

PRODUCT INFORMATION

(-)-3,4-Methylenedioxy Pyrovalerone (hydrochloride)

Item No. 14524

Formal Name: 1-(benzo[d][1,3]dioxol-5-yl)-2-(pyrrolidin-1-yl)pentan-1-one, monohydrochloride

Synonym: (-)-3,4-MDPV

MF: $C_{16}H_{21}NO_3 \cdot HCl$

FW: 311.8

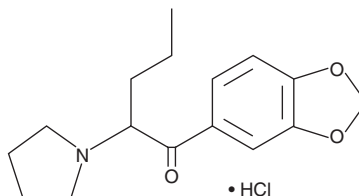
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 236, 283, 321 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

3,4-Methylenedioxy Pyrovalerone (3,4-MDPV) (Item No. 10624) is an analog of pyrovalerone (Item No. 10817) which includes the 3,4-methylenedioxy moiety found on 3,4-methylenedioxymethamphetamine, a DEA Schedule I controlled substance. This synthetic cathinone has been reported by the DEA to be abused as a central nervous system stimulant.^{1,2} It is an addictive stimulant that has distinct effects at different doses.³⁻⁵ (-)-3,4-MDPV is an enantiomerically pure isomer of 3,4-MDPV. Its properties, compared to the (+)-3,4-MDPV stereoisomer (Item No. 14525), have not been evaluated. (-)-3,4-Methylenedioxy Pyrovalerone is regulated as a Schedule 1 compound in the United States. This product is intended for forensic and research purposes.

References

1. Brandt, S.D., Freeman, S., Sumnall, H.R., *et al.* Analysis of NRG 'legal highs' in the UK: Identification and formation of novel cathinones. *Drug Test. Anal.* **3**(9), 569-575 (2011)
2. Prosser, J.M. and Nelson, L.S. The toxicology of bath salts: A review of synthetic cathinones. *J. Med. Toxicol.* **8**(1), 33-42 (2012).
3. Coppola, M. and Mondola, R. 3,4-Methylenedioxypyrovalerone (MDPV): Chemistry, pharmacology and toxicology of a new designer drug of abuse marketed online. *Toxicol. Lett.* **208**(1), 12-15 (2012).
4. Huang, P.-K., Aarde, S.M., Angrish, D., *et al.* Contrasting effects of d-methamphetamine, 3,4-methylenedioxymethamphetamine, 3,4-methylenedioxypyrovalerone, and 4-methylmethcathinone on wheel activity in rats. *Drug Alcohol Depend.* **126**(1-2), 168-175 (2012).
5. Watterson, L.R., Kufahl, P.R., Nemirovsky, N.E., *et al.* Potent rewarding and reinforcing effects of the synthetic cathinone 3,4-methylenedioxypyrovalerone (MDPV). *Addict. Biol.* **19**(2), 165-174 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/20/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM